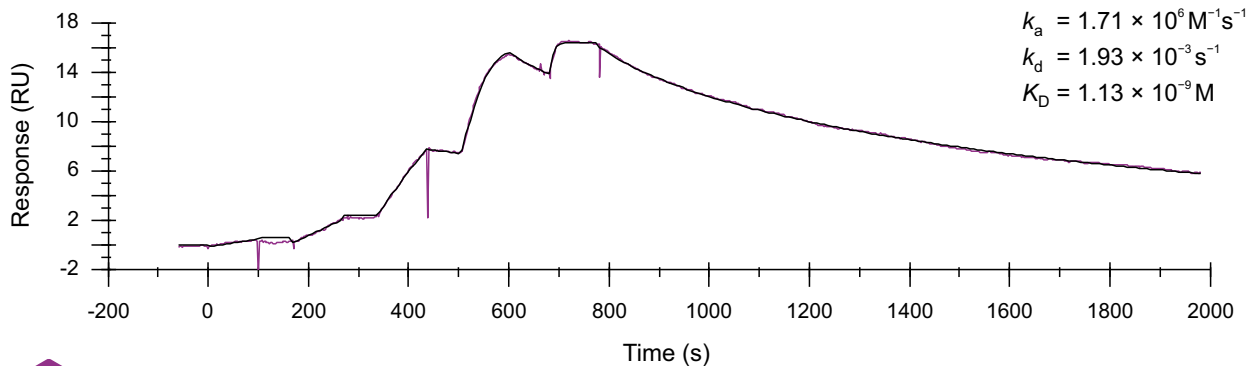


Binding kinetics of staurosporine on SYK



Oncolines

A SYMERES COMPANY

www.residencetimer.com
services@oncolines.com